

Rice seedlings OsMADS6 Primer pair

Cat. No. C17040009

Lot #: 001

Size: 50 µl/ 500 µl

Concentration: 10 µM

Specificity: Rice

Amplicon length: 140 base pairs

Format: 10 µM solution in MiliQ water (5 µM of each primer)

Storage: For long storage, store at -20°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: This primer pair specifically amplifies a genomic region from the MADS6 gene from Rice (*Oriza sativa* ssp Japonica cv. Nipponbare). The primers are thoroughly tested and optimized for routine SYBR® Green Real-Time qPCR assay following ChIP.

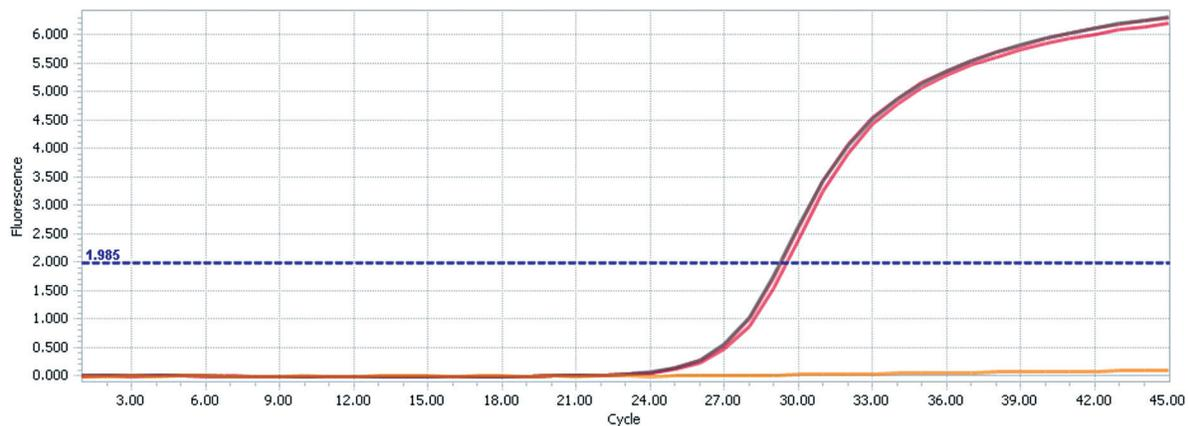


Figure 1.

Sheared DNA (0.5 ng) from rice seedlings (*Oriza sativa* ssp Japonica cv. Nipponbare) was analysed by real-time PCR using the Diagenode primers to amplify a genomic region in the MADS6 gene. 0.5 µl of provided primer pairs was used in a total volume of 10 µl. Real-time PCR was performed with the BioRad iCycler using SYBR Green. PCR conditions were as follows: an initial incubation at 95°C for 3 minutes, followed by 45 cycles of 30 seconds at 95°C, 30 seconds at 60°C and 30 seconds at 42°C. Triplicates are shown in red and grey. The dotted line represents the threshold position. The no template control is shown in orange and shows no amplification.

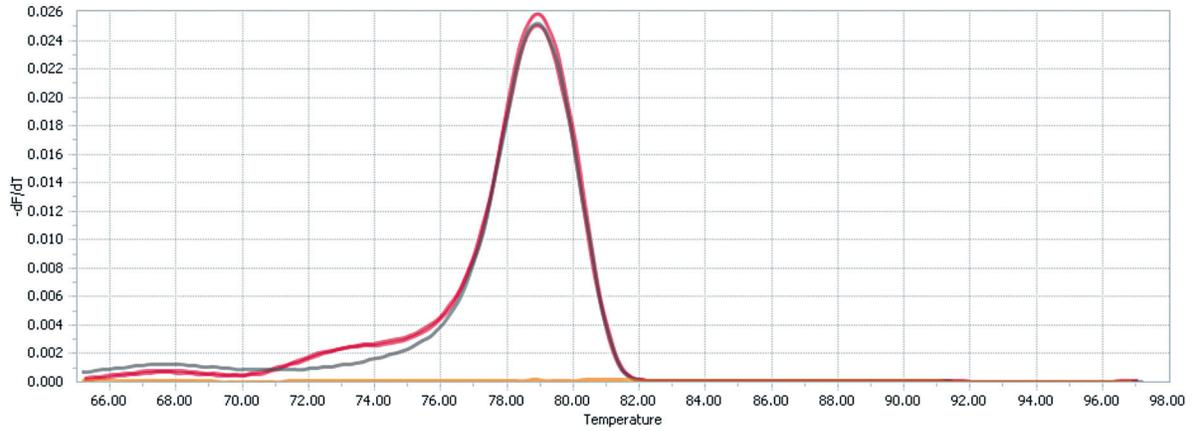


Figure 2.

Melting curve of PCR product amplified with the Rice seedlings OsMADS6 Primer pair (cat. No. C17040009). Real-time PCR was performed as described above. The melting curve analysis of the PCR product was performed by increasing the temperature from 55°C to 95°C in 0.5°C increments. No amplification was found in the negative control (orange).

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
 Rue Bois Saint-Jean, 3
 4102 Seraing (Ougrée) - Belgium
 Tel: +32 4 364 20 50
 Fax: +32 4 364 20 51
 orders@diagenode.com
 info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
 Denville, NJ 07834 - USA
 Tel: +1 862 209-4680
 Fax: +1 862 209-4681
 orders.na@diagenode.com
 info.na@diagenode.com

Last update: February 17, 2016